

Rabbit Anti-Alexa Fluor 647 Antibody, iFluor® 488-Labeled

Catalog Number: 46565

Unit Size: 100 ug

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, BSA

Unit Details

Reconstitution Volume	0.1 mL ddH2O
-----------------------	--------------

Antibody Properties

Clonality	Polyclonal
Host	Rabbit
Immunogen	Alexa Fluor 647 coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 647 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 488 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 488
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

Rabbit Anti-Alexa Fluor 647 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules and is labeled with iFluor® 488, a green fluorescent dye with excitation/emission of ~491/516 nm. This direct labeling format allows one-step detection of Alexa Fluor® 647-labeled targets without the need for secondary antibodies. The antibody may cross-react with related fluorophores such as iFluor® 647, Cy5, ATTO 647, and DyLight® 650. Its green fluorescence is spectrally distinct from far-red targets, making it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection. *Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.*